

Independent Auditors' Fees

Principal Independent Auditors

KPMG S.A.

480, avenue du Prado – CS 90021 – 13272 Marseille Cedex 08 – France

Represented by John Evans.

First commissioned by the Annual General Shareholders' Meeting on 7 April 2015.

Date commission expires: 2021 Annual General Shareholders' Meeting to approve the 2020 financial statements.

Member of the Compagnie régionale de Versailles.

Deloitte et Associés

7, boulevard Jacques Saadé - Quai de la Joliette – 13235 Marseille Cedex 2 – France

Represented by Philippe Battisti.

First commissioned by the Annual General Shareholders' Meeting on 19 May 2006.

Date commission expires: 2024 Annual General Shareholders' Meeting to approve the 2023 financial statements.

Member of the Compagnie régionale de Versailles.

Independent Auditors' Fees

€ in K	2019		KPMG 2018		2019		Deloitte 2018	
Audit								
Independent audit, certification, parent company & consolidated financial statements								
Parent company	63	8.3%	61	7.4%	55	36.9%	54	37.2%
Subsidiaries	799	90.4%	754	92.0%	94	63.1%	91	62.8%
Services directly related to audit services								
Parent company								
Subsidiaries								
Subtotal	862	98.6%	815	99.5%	149	100.0%	145	100.0%
Other services								
Legal, tax, corporate	12	1.4%	5	0.5%	0	0.0%	0	0.0%
Information technology, other	0	0.0%	0	0.0%	0	0.0%	0	0.0%
Subtotal	12	1.4%	5	0.5%	0	0.0%	0	0.0%
Total	874	100.0%	820	100%	149	100.0%	145	100%

Substitute Independent Auditors

SALUSTRO REYDEL

Tour Eqho -2 avenue Gambetta -92066 Paris La
Défense Cedex

First commissioned by the Annual General Share-
holders' Meeting on 7 April 2015.

Date commission expires: 2021 Annual General
Shareholders' Meeting to approve the 2020 financial
statements.

Member of the Compagnie régionale de Versailles.

	2019	Other		2019		Total	
		2018				2018	
				118	9.5%	115	10.2%
76	35.0%	60	37.8%	969	78.1%	905	80.6%
76	35.0%	60	37.8%	1,087	87.6%	1,020	90.8%
117	53.5%	51	32.4%	129	10.4%	56	5.0%
25	11.5%	47	29.7%	25	2.0%	47	4.2%
142	65.0%	98	62.2%	154	12.4%	103	9.2%
218	100.0%	158	100%	1,241	100.0%	1,123	100%